

## HIPK4

(Homeodomain-interacting protein kinase 4)

- **CAT#:** HIPK4
- **Enzyme:** Human HIPK4
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	621
Ro 31-8220	961
GW 5074	882
Rottlerin	61,400

- **Scientific Information:** Q8NE63 (<http://www.expasy.org/uniprot/Q8NE63>)

### Reference compound IC50 for HIPK4

